

ALDH1A2 Antibody [#C18344]

Catalog Number:	C18344
Concentration:	1mg/ml
Swiss-Prot No.:	O94788
Other Names:	AL1A2; aldehyde dehydrogenase 1 family, member A2; aldehyde dehydrogenase 1A2; aldehyde dehydrogenase family 1 member A2; EC 1.2.1.36; II; RALDH; RALDH 2; RALDH(II); RALDH2; RALDH2-T; retinal dehydrogenase 2; retinaldehyde dehydrogenase 2; retinaldehyde-specific dehydrogenase type 2
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg ²⁺ and Ca ²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized peptide derived from internal of human ALDH1A2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ALDH1A2 antibody detects endogenous levels of total ALDH1A2 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Ono Y., Mol. Cell. Biol. 18:6939-6950(1998). Livingston R.J., Submitted (DEC-2005) to the EMBL/GenBank/DDBJ databases. The MGC Project Team; Genome Res. 14:2121-2127(2004).